

Certificate of Analysis

MSK2, active

(Recombinant enzyme expressed in Sf21 insect cells)

Item # 14-616, 14-616-K, 14-616M

Parent Lot # D7PN035U

The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialling runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialling run.

Product Description: N-terminal 6His-tagged, recombinant, human MSK2, amino acids 2–end, expressed by baculovirus in Sf21 insect cells. Purified using Ni²⁺/NTA-agarose. Activated using MAPK2 (cat# 14-173) and repurified using Ni²⁺/NTA-agarose. Purity 64% by SDS-PAGE and Coomassie blue staining. MW = 89.9kDa.

Specific Activity (Parent lot# D7PN035U): 130U/mg, where one unit of MSK2, active activity is defined as 1nmol phosphate incorporated into 30µM modified Crosstide (GRPRTSSFAEGKK) per minute at 30°C with a final ATP concentration of 100µM.

Formulation: 0.9285mg/ml of enzyme in 50mM Tris/HCl pH7.5, 300mM NaCl, 0.1mM EGTA, 1mM benzamidine, 0.2mM PMSF, 0.1% 2-mercaptoethanol, 0.03% Brij-35, 270mM sucrose. Frozen solution.

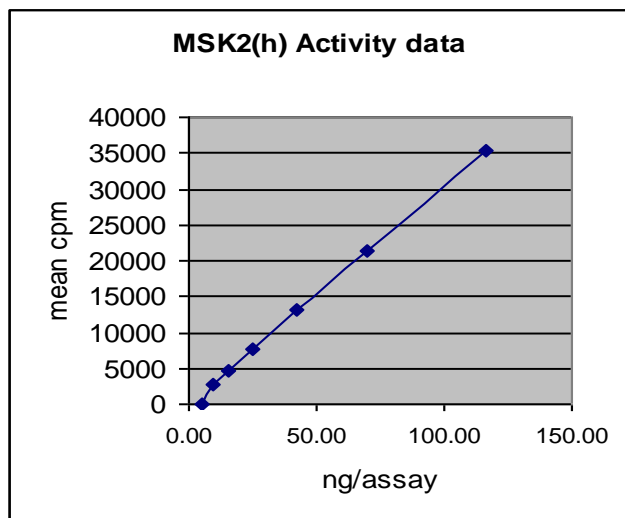
Storage and Stability: On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

Handling Recommendations: Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled micro-centrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

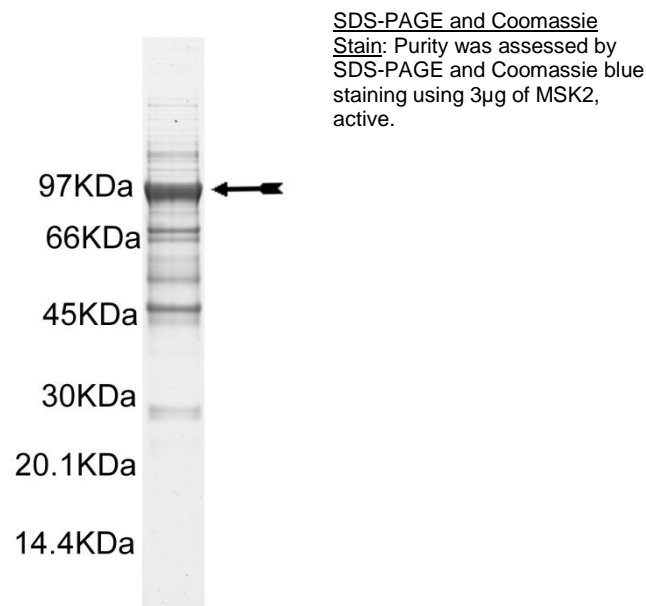
**FOR IN VITRO RESEARCH USE ONLY
NOT FOR USE IN HUMANS OR ANIMALS**

Quality Control Testing

Kinase Assay: 42–116ng of this lot of enzyme phosphorylated 30µM modified Crosstide (GRPRTSSFAEGKK) in the assay described on page two. Assay background was subtracted from the actual counts to yield the results shown below.



MS Tryptic Fingerprint: Confirmed identity as MSK2 with the translated sequence listed on page three.



Certificate of Analysis

Kinase Assay Protocol

Stock Solutions:

1. **5 x Reaction Buffer:** 40mM MOPS/NaOH pH7.0, 1mM EDTA.
2. **Modified Crosstide (GRPRTSSFAEGKK):** Use a final assay concentration of 30 μ M. Make a 300 μ M stock. Add 2.5 μ l of stock per assay point.
3. **MSK2, active:** Dilute with 20mM MOPS/NaOH pH7.0, 1mM EDTA, 5% glycerol, 0.01% Brij-35, 0.1% 2-mercaptoethanol, 1mg/ml BSA. Use 42–116ng per assay point.
4. **[γ -³³P]ATP:** 2.5x magnesium acetate/[γ -³³P]ATP cocktail: 25mM MgAc and 0.25mM ATP to which is added [γ -³³P]ATP (specific activity approximately 500 - 800cpm/pmol as required.)

Assay Procedure (96 well plate format):

1. Add 5 μ l of 5 x reaction buffer per assay to wells.
2. Add 2.5 μ l of **Crosstide (GRPRTSSFAEGKK)**.
3. Add **2.5 μ l (42–116ng) MSK2, active**.
4. Add 5 μ l of dH₂O.
5. Add 10 μ l of diluted [γ -³³P]ATP mixture.
6. Incubate for 10 minutes at 30°C.
7. Stop the reaction by adding 5 μ l of 3% phosphoric acid.
8. Transfer a 10 μ l aliquot onto the appropriate area of a **P30 Filtermat**.
9. Wash the filtermat three times for 5 minutes with 75mM phosphoric acid.
10. Wash the filtermat once for 2 minutes with methanol.
11. Transfer the filtermat to a sealable plastic bag and add 4ml of scintillation cocktail.
12. Read in a scintillation counter. Compare cpm of enzyme samples with cpm of control samples that contain all assay components plus 1 μ l of 30% phosphoric acid.

Certificate of Analysis

MSK2 Sequence Information

Protein	Human MSK2
Tags	N-terminal 6His
Native sequence	G37 of the recombinant protein is equivalent to G2 of human MSK2
Accession number	GenBank AJ010119

Recombinant MSK2 amino acid sequence:

```

1 MSYYHHHHHH DYDIPTTENL YFQGAMGIRN SKAYVDGDED DDESCAVELR ITEANLTGHE
61 EKVSVENFEL LKVLGTGAYG KVFLVRKAGG HDAGKLYAMK VLRKAALVQR AKTQEHTRTE
121 RSVLELVRQA PFLVTLHYAF QTDAKLHLIL DYVSGGEMFT HLYQRQYFKE AEVRVYGGEI
181 VLALEHLHLK GIIYRDLKLE NVLLDSEGI VLTDFGLSKE FLTEEKERTF SFCGTIEYMA
241 PEIIRSKTGH GKAVDWWSLG ILLFELLTGA SPFTLEGERN TQAEVSRRIL KCSPPFPRI
301 GPVAQDLLQR LLCKDPKKRL GAGPQGAQEV RNHPFFQGLD WVALAARKIP APFRPQIRSE
361 LDVGNFAEEF TRLEPVYSPP GSPPPDPRI FQGYSFVAPS ILFDHNNAVM TDGLEAPGAG
421 DRPGRAAVAR SAMMQDSPFF QQYELDLREP ALGQGSFVVC RRCRQRQSGQ EFAVKILSRR
481 LEANTQREVA ALRLCQSHPN VVNLHEVHHD QLHTYLVLEL LRGGELLEHI RKKRHFSESE
541 ASQILRSLVS AVSFMHEEAG VVHRDLKPEN ILYADDTPGA PVKIIDFGFA RLRPQSPGVP
601 MQTPCFTLQY AAPELLAQQG YDESCDLWSL GVILYMMLSG QVPFQGASGQ GGQSQAAEIM
661 CKIREGRFSL DGEAWQGVSE EAKELVRGLL TVDPAKRLKL EGLRGSSWLQ DGSARSSPPL
721 RTPDVLESSG PAVRSGLNAT FMAFNRGKRE GFFLKSVENA PLAKRRKQKL RSATASRRGS
781 PAPANPGRAP VASKGAPRA NGPLPPS

```

Recombinant MSK2 nucleotide sequence:

```

1 atgtcgtact accatcacca tcaccatcac gattacgata tcccaacgac cgaaaacctg
61 tattttcagg gcgccatggg gatccggaat tcaaaggcct acgtcgacgg ggacgaggac
121 gacgatgaga gctgcgccgt ggagctgcgg atcaccgaag ccaacctgac cgggcacgag
181 gagaaggtga gcgtaggaga cttcgagctg ctcaaggtgc tgggcacggg agcctacggc
241 aaggtgttcc tggtagcggaa ggcgggcggg cacgacgagg ggaagctgta cgccatgaag
301 gtgctgcgca aggcagcgcct ggtgcagcgc gccaaagacg aagagcacac ggcgaccgag
361 cgctcggtgc tggagctggt gcgccaggcg ccttctctgg tcacgctgca ctacgctttc
421 cagacggatg ccaagctgca cctcatcctg gactatgtga gcggcgggga gatgttcacc
481 cacctctacc agcgccagta cttcaaggag gctgaggtgc gcgtgatagg gggtagatc
541 gtgctggccc tggaaacact gcacaagctc ggcacattt accgagacct gaaactggag
601 aatgtgctgc tggactccga gggccacatt gcctctacgg acttcgggct gagcaaggag
661 ttctgtagcg aggagaaaga gcggaccttc tccttctgtg gcaccatcga gtacatggcc
721 cccgaaatca tccgtagcaa gacggggcat ggcaaggctg tggactggtg gaggcctggc
781 atcttgcctc tcgagctgct gacgggggccc tcgcccttca ccctggaggg cgagaggaac
841 acgcaggctg aggtgtctcg acggatcctg aagtgtctcc ctcccttccc ccctcggatc
901 gggcccgtgg cgcaggacct gctgcagcgg ctgctttgta aggatcctaa gaagcgattg
961 ggcgcggggc cccagggggc acaagaagtc cggaaccatc ctttcttcca gggcctcgat
1021 tgggtggctc tggctgccag gaagattcca gccccattcc ggccccaaat ccgctcagag
1081 ctggatgtgg gcaactttgc ggaggaattc actcggctgg agcctgtcta ctaccccct
1141 ggcagcccc cacctgggga cccccgaatc tttcagggat actcctttgt ggcaccctcc
1201 attctctttg accacaacaa cgcggtgatg accgatgggc tggaaagcgc tgggtgctgga
1261 gaccggccag gtcgggcagc ggtggccagg agcgctatga tgcaggactc gcccttcttc
1321 cagcagtacg agctggacct gcgggagcct gcgctgggccc agggcagctt ttctgtgtgt
1381 cgccgctgcc gccagcgcca gageggccag gagttcgag tcaagatcct cagtcgcagg
1441 ctggaggcga acacgcagcg cgaagtggct gccctgcgcc tgtgccagtc acacccaac
1501 gtggtgaatc tgcacgaggt gcatcacgac cagctgcaca cgtacctggt cctggagctg
1561 ctgcggggcg gggagctgct ggagcacatc cgcaagaagc ggcacttcag cgagtcggaa
1621 gcaagccaga tcctgcgcag cctcgtgtcg gccgtgagct tcatgcacga ggaggcgggc

```

Certificate of Analysis

```
1681 gtggtgcacc gcgacctcaa gccggagaac atcctgtacg ccgacgacac gcccggggcc
1741 ccggtgaaaa tcatcgactt cgggttcgcg cggttgcggc cgcagagtcc cgggggtgcc
1801 atgcagacgc cctgcttcac gctgcagtac gctgcccccg agctgctggc gcagcagggc
1861 tacgacgagt cctgcgacct ctggagcctg ggcgtcattc tgtacatgat gctgtcgggg
1921 caggtcccct tccagggggc ctctggccag ggcgggcaga gccaggcggc cgagatcatg
1981 tgcaaaatcc gcgagggggc cttctccctt gacggggagg cctggcaggg tgtatccgag
2041 gaagccaagg agctggtccg agggctcctg accgtggacc ccgccaagcg gctgaagctc
2101 gagggactgc ggggcagctc gtggctgcag gacggcagcg cgcgctcctc gccccgctc
2161 cggacgcccc acgtgctcga gtctctggtg cccgcagtgc gctcgggtct caacgccacc
2221 ttcattggcat tcaaccgggg caagcgggag ggcttcttcc tgaagagcgt ggagaatgca
2281 cccctggcca agcggcggaa gcagaagctg cggagcgcca ccgcctccc cgggggctcc
2341 cctgcaccag ccaaccggg ccgagcccc gtcgcctcca aaggggcccc ccgccgagcc
2401 aacggcccc tgccccctc ctaa
```

Reviewed and approved by site quality representative.

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

© 2014 Eurofins Pharma Discovery Services UK Limited is an independent member of Eurofins Discovery Services